

Helenalin

NF- κ B inhibitor

Potent anti-inflammatory and antineoplastic agent. Induces a CD95 death receptor-independent apoptosis even when the mitochondria protecting proteins Bcl-X_L or Bcl-2 are overexpressed. Also an inhibitor of NF- κ B.

Citations: 8

[View Online »](#)

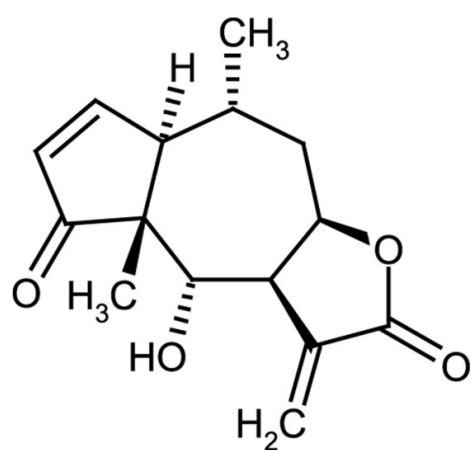
Ordering Information

[Order Online »](#)

ALX-350-120-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	6754-13-8
Couple Target	NF-kappaB
Couple Type	Inhibitor
Formula	$C_{15}H_{18}O_4$
Identity	Determined by NMR.
MW	262.3
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol.
Source	Isolated from <i>Arnica</i> species.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)